

# PRODUCT INFORMATION



## Z-VDVAD-FMK

Item No. 17484

CAS Registry No.: 210344-92-6

Formal Name: N-[(phenylmethoxy)carbonyl]-L-valyl-L- $\alpha$ -aspartyl-L-valyl-N-[(1S)-3-fluoro-1-(2-methoxy-2-oxoethyl)-2-oxopropyl]-L-alaninamide, methyl ester

Synonyms: Z-Val-Asp-(OMe)-Val-Ala-Asp-(OMe)-FMK, Z-VDVAD-Fluoromethyl Ketone

Peptide Sequence: Z-VXVAX-FMK (X = O-methylated aspartic acid)

MF:  $C_{32}H_{46}FN_5O_{11}$

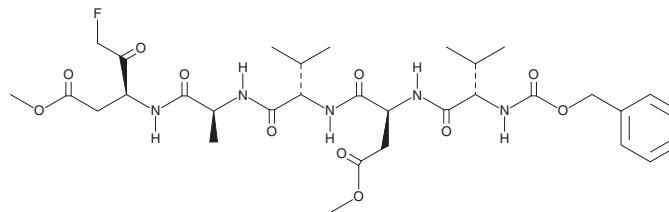
FW: 695.7

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Z-VDVAD-FMK is supplied as a solid. A stock solution may be made by dissolving the Z-VDVAD-FMK in the solvent of choice, which should be purged with an inert gas. Z-VDVAD-FMK is soluble in the organic solvent DMSO.

### Description

Z-VDVAD-FMK is a peptide caspase-2 inhibitor.<sup>1,2</sup> Intravitreal injection of Z-VDVAD-FMK (4,000 ng/ml) reduces optic nerve crushing-induced retinal ganglion cell apoptosis in rats.<sup>1</sup> It also prevents endothelial cell apoptosis in a dog model of subarachnoid hemorrhage-induced chronic cerebral vasospasm.<sup>2</sup>

### References

1. Vigneswara, V., Berry, M., Logan, A., *et al.* Pharmacological inhibition of caspase-2 protects axotomised retinal ganglion cells from apoptosis in adult rats. *PLoS One* **7(12)**, e53473 (2012).
2. Aoki, K., Zubkov, A.Y., Ross, I.B., *et al.* Therapeutic effect of caspase inhibitors in the prevention of apoptosis and reversal of chronic cerebral vasospasm. *J. Clin. Neurosci.* **9(6)**, 672-677 (2002).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/04/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM